



# Filtration of wine samples and measure of the wine clogging index

## Product Information

When small volumes of wine need to be filtered for their presentation and tasting, MidiCaps<sup>®</sup> capsules simplify your filtration in complete safety and while respecting your wines.



### From a few liters to hundreds of liters!

The filter materials of the MidiCaps<sup>®</sup> are the same as those used on the bottling lines but with surface areas 9 times smaller.

### Applications

According to the filter materials chosen and their pore size:

- Clarification
- Particle retention
- Bioburden reduction
- Sterile filtration

MidiCaps<sup>®</sup> filtration capsules simplify your filtration in complete safety and while respecting your wines.

### MidiCaps<sup>®</sup>: ready-to-use filtration capsules

#### Traceability

Each capsule is identified: type of filter, reference, pore size and batch number.

Engraving on the capsules of the conditions for use and direction of filtration.

#### Quality

Purge and sanitary type vent to optimize the use of the capsules. MidiCaps<sup>®</sup> have a validation guide according to FDA recommendations. Quality assurance certificates in each box.

#### Ease of use

Available with different connections: triclamp or hose barb for flexible tubing.

Sterilized by autoclaving or hot water. Contact us for advice on the regeneration and storage of capsules.

## Filtration of wine samples and measurement of the clogging index of wine



### 5 liter vessel

Sartorius provides a stainless steel vessel for the filtration of small volumes of wine.

### Application

- Filtration of wine samples with filtration capsules
- Measurement of the clogging index and Vmax

### Practical and versatile

#### Ease of use

- Stainless steel system mounted on an height-adjustable scaffold
- 0 to 6 bar pressure gauge.
- Two valves at the highest point for the intake of air and release into the atmosphere after filtration.
- A valve at the lowest point for the outlet of wine.



### 2 in 1!

- Equipped as standard with a support disc of 25 mm in diameter to measure the clogging index and Vmax.
- Triclamp connector for mounting a capsule filter and filtering wine samples.

## Technical Specifications

Dimensions (W × D × H )	450 × 460 × 1300 mm (17 3/4 × 18 1/8 × 51 3/16 in)
Weight (approx)	17 kg (37.48 lbs)

## Ordering Information

### Vessel

Description	Reference
Complete 5 liter vessel	FR-INDICE5L349

### Membrane for the measure of the clogging index and Vmax

Description	Reference
0.65 µm cellulose acetate membrane 25 mm in diameter (box of 100)	111 05 025N

### MidiCaps® capsules for filtration of samples

Filter material	Applications	Pore size	Inlet connection	Inlet connection	References
Positively charged fiberglass	Brightness of the wines	0.65 µm	Clamp	Hose barb	5555305P9--SO--A
			Hose barb	Hose barb	5555305P9--00--A
		1.2 µm	Clamp	Hose barb	5555303P9--SO--A
			Hose barb	Hose barb	5555303P9--00--A
Polypropylene	Clarification of the wines and particle retention	0.65 µm	Clamp	Hose barb	5055305P9--SO--A
			Hose barb	Hose barb	5055305P9--00--A
		3 µm	Clamp	Hose barb	5055302P9--SO--A
			Hose barb	Hose barb	5055302P9--00--A
			Clamp	Clamp	5055302P9--SS--A
		5 µm	Hose barb	Hose barb	5055342P9--00--A
Cellulose acetate	Sterilizing filtration	0.65 µm	Clamp	Clamp	5625305A9--SS--A
		0.45 µm	Clamp	Hose barb	5625306G9--SO--A

# Sales and Service Contacts

For further contacts, visit [www.sartorius-stedim.com](http://www.sartorius-stedim.com)

## Europe

### Germany

Sartorius Stedim Biotech GmbH  
August-Spindler-Strasse 11  
37079 Goettingen

Phone +49.551.308.0

Sartorius Stedim Systems GmbH  
Robert-Bosch-Strasse 5-7  
34302 Guxhagen

Phone +49.5665.407.0

### France

Sartorius Stedim FMT S.A.S.  
ZI des Paluds  
Avenue de Jouques - CS 91051  
13781 Aubagne Cedex

Phone +33.442.845600

Sartorius Stedim France SAS  
ZI des Paluds  
Avenue de Jouques - CS 71058  
13781 Aubagne Cedex

Phone +33.442.845600

### Austria

Sartorius Stedim Austria GmbH  
Modectcenterstrasse 22  
1030 Vienna

Phone +43.1.7965763.18

### Belgium

Sartorius Stedim Belgium N.V.  
Rue Colonel Bourg 105  
1030 Bruxelles

Phone +32.2.756.06.80

### Hungary

Sartorius Stedim Hungária Kft.  
Kagyló u. 5  
2092 Budakeszi

Phone +36.23.457.227

### Italy

Sartorius Stedim Italy S.r.l.  
Via dell'Antella, 76/A  
50012 Antella-Bagno a Ripoli (FI)

Phone +39.055.63.40.41

### Netherlands

Sartorius Stedim Netherlands B.V.

Phone +31.30.60.25.080

[filtratie.nederland@sartorius-stedim.com](mailto:filtratie.nederland@sartorius-stedim.com)

### Poland

Sartorius Stedim Poland Sp. z o.o.  
ul. Wrzesinska 70  
62-025 Kostrzyn

Phone +48.61.647.38.40

### Russian Federation

LLC "Sartorius Stedim RUS"  
Vasilyevsky Island  
5<sup>th</sup> line 70, Lit. A  
199178 St. Petersburg

Phone +7.812.327.53.27

### Spain

Sartorius Stedim Spain, S.A.U.  
Avda. de la Industria, 32  
Edificio PAYMA  
28108 Alcobendas (Madrid)

Phone +34.913.586.098

### Switzerland

Sartorius Stedim Switzerland AG  
Ringstrasse 24 a  
8317 Tagelswangen

Phone +41.52.354.36.36

### U.K.

Sartorius Stedim UK Ltd.  
Longmead Business Centre  
Blenheim Road, Epsom  
Surrey KT19 9 QQ

Phone +44.1372.737159

### Ukraine

LLC "Sartorius Stedim RUS"  
Post Box 440 "B"  
01001 Kiev, Ukraine

Phone +380.44.411.4918

## Africa

### Republic of South Africa

Sartorius South Africa (Pty) Ltd  
Unit 4, Alphen Square South  
853 16th Road  
Midrand 1685, RSA

Phone +27.11.315.5444

## Americas

### USA

Sartorius Stedim North America Inc.  
5 Orville Drive, Suite 200  
Bohemia, NY 11716

Toll-Free +1.800.368.7178

### Argentina

Sartorius Argentina S.A.  
Int. A. Ávalos 4251  
B1605ECS Munro  
Buenos Aires

Phone +54.11.4721.0505

### Brazil

Sartorius do Brasil Ltda  
Avenida Senador Vergueiro 2962  
São Bernardo do Campo  
CEP 09600-000 - SP- Brasil

Phone +55.11.4362.8900

### Mexico

Sartorius de México, S.A. de C.V.  
Libramiento Norte de Tepotzotlan s/n,  
Colonia Barrio Tlacateco,  
Municipio de Tepotzotlan,  
Estado de México,  
C.P. 54605

Phone +52.55.5562.1102

[leadsmex@sartorius.com](mailto:leadsmex@sartorius.com)

## Asia | Pacific

### Australia

Sartorius Stedim Australia Pty. Ltd.  
Unit 5, 7-11 Rodeo Drive  
Dandenong South Vic 3175

Phone +61.3.8762.1800

### China

Sartorius Stedim (Shanghai)  
Trading Co., Ltd.  
3rd Floor, North Wing, Tower 1  
No. 4560 Jinke Road  
Zhangjiang Hi-Tech Park  
Pudong District  
Shanghai 201210, P.R. China

Phone +86.21.6878.2300

Sartorius Stedim (Shanghai)  
Trading Co., Ltd.  
Beijing Branch Office  
No. 33 Yu'an Road  
Airport Industrial Park Zone B  
Shunyi District, Beijing 101300

Phone +86.10.8042.6501

Sartorius Stedim (Shanghai)  
Trading Co., Ltd.  
Guangzhou Branch Office  
Room 1105  
Xing Guang Ying Jing Building  
No. 119, Shui Yin Road  
Yue Xiu District, Guangzhou 510075

Phone +86.20.3836.4193

### India

Sartorius Stedim India Pvt. Ltd.  
#69/2-69/3, NH 48, Jakkasandra  
Nelamangala Tq  
562 123 Bangalore, India

Phone +91.80.4350.5250

### Japan

Sartorius Stedim Japan K.K.  
4th Fl., Daiwa Shinagawa North Bldg.  
8-11, Kita-Shinagawa 1-chome  
Shinagawa-ku, Tokyo, 140-0001 Japan

Phone +81.3.4331.4300

### Malaysia

Sartorius Stedim Malaysia Sdn. Bhd.  
Lot L3-E-3B, Enterprise 4  
Technology Park Malaysia  
Bukit Jalil  
57000 Kuala Lumpur, Malaysia

Phone +60.3.8996.0622

### Singapore

Sartorius Stedim Singapore Pte. Ltd.  
10 Science Park Rd  
The Alpha #02-13/14  
Singapore Science Park II  
Singapore 117684

Phone +65.6872.3966

### South Korea

Sartorius Korea Biotech Co., Ltd.  
8th Floor, Solid Space B/D,  
PanGyoYeok-Ro 220, BunDang-Gu  
SeongNam-Si, GyeongGi-Do, 463-400

Phone +82.31.622.5700



▶ [www.sartorius-stedim.com](http://www.sartorius-stedim.com)